

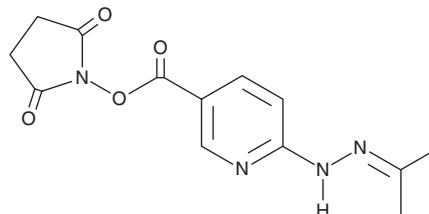
PRODUCT INFORMATION



HyNic NHS ester

Item No. 45697

CAS Registry No.: 362522-50-7
Formal Name: 6-[2-(1-methylethylidene)hydrazinyl]-3-pyridinecarboxylic acid, 2,5-dioxo-1-pyrrolidinyl ester
Synonyms: HyNic N-hydroxysuccinimide ester, S-HyNic, SANH
MF: C₁₃H₁₄N₄O₄
FW: 290.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

HyNic NHS ester is supplied as a solid. A stock solution may be made by dissolving the HyNic NHS ester in the solvent of choice, which should be purged with an inert gas. HyNic NHS ester is sparingly soluble (1-10 mg/ml) in DMSO.

Description

HyNic NHS ester is a heterobifunctional linker molecule.¹ It contains an amine-reactive N-hydroxysuccinimide (NHS) moiety and a hydrazine moiety that reacts with the aldehyde functional group in SFB (succinimidyl 4-formylbenzoate; Item No. 30477) to form a *bis*-aryl hydrazone linker that is quantifiable spectrophotometrically at 354 nm. HyNic NHS ester has been used in the formation of small molecule-antibody, polymer-enzyme, protein-liposome, and peptide-peptide conjugates.¹⁻⁴

References

1. Grotzky, A., Manaka, Y., Kojima, T., *et al.* Preparation of catalytically active, covalent α-polylysine-enzyme conjugates via UV/vis-quantifiable bis-aryl hydrazone bond formation. *Biomacromolecules* **12**(1), 134-44 (2011).
2. Gardner, M.W. and Brodbelt, J.S. Ultraviolet photodissociation mass spectrometry of bis-aryl hydrazone conjugated peptides. *Anal. Chem.* **81**(12), 4864-4872 (2009).
3. Gadd, J.R., Greco, F., Cobb, A.J.A., *et al.* Targeted activation of toll-like receptors: Conjugation of a toll-like receptor 7 agonist to a monoclonal antibody maintains antigen binding and specificity. *Bioconjug. Chem.* **26**(8), 1743-1752 (2015).
4. Nagata, H., Yoshimoto, M., and Walde, P. Preparation and catalytic properties of carbonic anhydrase conjugated to liposomes through a bis-aryl hydrazone bond. *ACS Omega* **8**(21), 18637-17652 (2023).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/26/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM